

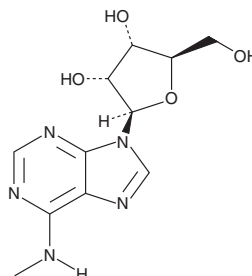
# PRODUCT INFORMATION



## N<sup>6</sup>-Methyladenosine

Item No. 28833

**CAS Registry No.:** 1867-73-8  
**Formal Name:** N-methyl-adenosine  
**Synonym:** NSC 29409  
**MF:** C<sub>11</sub>H<sub>15</sub>N<sub>5</sub>O<sub>4</sub>  
**FW:** 281.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 266 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

N<sup>6</sup>-Methyladenosine is an adenosine analog.<sup>1</sup> It inhibits epinephrine-induced contraction of isolated guinea pig ileum and thoracic aorta when used at a concentration of 10 μM. N<sup>6</sup>-Methyladenosine (1 mg/kg, i.v.) decreases arterial blood pressure and renal blood flow and increases peripheral resistance in anesthetized dogs.<sup>2</sup> It also inhibits tumor growth in the C3H/ST and C3HB/ST mouse models of spontaneous mammary adenocarcinomas.<sup>3</sup> N<sup>6</sup>-Methyladenosine is also the most prevalent mRNA modification in eukaryotes and has roles in cell viability and development.<sup>4</sup>

### References

1. Leslie, S.W., Borowitz, J.L., and Miya, T.S. Adenosine analogs: Structure-activity relationships in vascular and intestinal smooth muscle. *J. Pharm. Sci.* **62**(9), 1449-1452 (1973).
2. Bhanalaph, T., Mittelman, A., Ambrus, J.L., et al. Effect of adenosine and some of its analogs on renal hemodynamics. *J. Med.* **4**(3), 178-188 (1973).
3. Strong, L.C. and Matsunaga, H. Comparison of the effect of three nucleosides and a liver emulsion upon the inhibition of cancer in mice. *J. Surg. Oncol.* **4**(3), 248-254 (1972).
4. Wang, X., Lu, Z., Gomez, A., et al. N<sup>6</sup>-Methyladenosine-dependent regulation of messenger RNA stability. *Nature* **505**(7481), 117-120 (2014).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/04/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)